

Product Number:	103186
Lot Number:	502305
Expiration:	17-Jan-2021
Platform(s):	HD-1/HD-x, SRx

Component	Part Number	Lot Number
Bead Reagent	102246	006304
Detector Reagent	102248	006206
SBG Reagent	102250	006205
Sample Diluent	102252	006310
Calibrators	102255	933601
Controls	102256 - 102257	933901
RGP Reagent	103159	N/A ¹

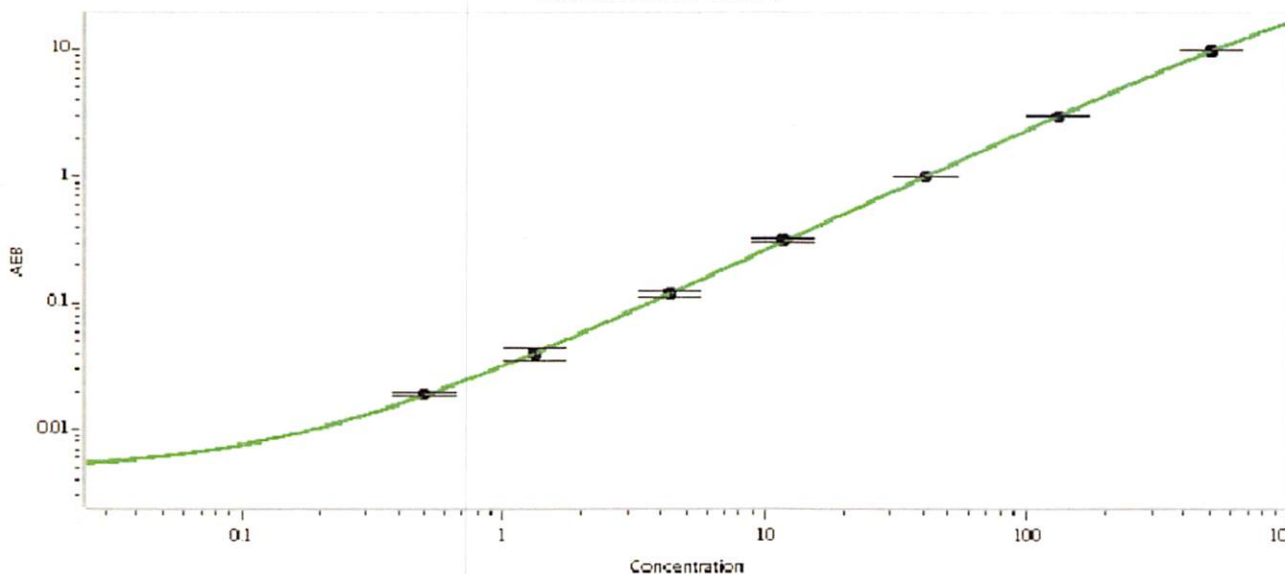
¹ RGP is not Kit Lot Specific



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*				
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	933901	3.29	2.56-4.80	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Control 2	933901	157	121-181	
Panel 1	922601	17.9	14.1-21.1	
Panel 2	922602	6.08	4.94-7.41	
Panel 3	922603	38.6	28.4-42.6	
Panel 4	922604	173	129-204	

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.501
C	1.32
D	4.32
E	11.8
F	41.1
G	131
H	506

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	26 Nov 2020
	Name	Title	Signature/Date	